

Live-Dead cell staining kit

Kit utilizes Live-dye™, a cell permeable green fluorescent dye (Ex(max): 488nm; Em(max): 518nm), to stain live cells. Dead cells can be easily stained by propidium iodide (PI), a red fluorescent dye (Ex(max): 488nm; Em(max): 615nm), which in viable cells is actively pumped out of the cytoplasm. Stained cells can be visualized by fluorescence microscopy using a band-pass filter (detects FITC and rhodamine). Cyto-dye has also been used to discriminate between apoptotic and non-apoptotic cells. *Ready-to-use.*

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-249-KI01	1Kit
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from light.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescence microscopy, Fluorescent detection

Quantity 100 tests



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)